


TG GRA7/p29 (residues 24-100)

Toxoplasma Gondii Dense Granule Protein 7
Toxoplasma gondii, recombinant, *E. coli*

Cat. No.	Amount
PR-1245	100 µg

For general laboratory use.

Shipping: shipped on gel packs

Storage Conditions: store at -20 °C

Additional Storage Conditions: avoid freeze/thaw cycles

Shelf Life: 12 months

Purity: > 95 % (SDS-PAGE)

Form: liquid (Supplied in 50 mM Tris-HCl pH 8.0, 1.5 M urea and 50% glycerol)

Applications:

Antigen in ELISA and Western blots, excellent antigen for detection of *Toxoplasma gondii* with minimal specificity problems.

Description:

The protein contains the p29 (GRA7) immunodominant regions, amino acids 24-100. The protein is purified by proprietary chromatographic technique.

Background: *Toxoplasma gondii* is an obligate intracellular protozoan parasite that infects all warmblooded animals, including humans, and causes toxoplasmosis. GRA7 is a recently discovered 29-kDa protein. Like GRA1, it is secreted from the dense granules, and it reacts with sera from humans with acute and chronic infections.

Specificity: Immunoreactive with sera of *T. gondii*-infected individuals.

Selected References:

Hiszczynska-Sawicka *et al.* (2003) High yield expression and singlestep purification of *Toxoplasma gondii* SAG1, GRA1, and GRA7 antigens in *Escherichia coli*. *Protein Expr. Purif.* **27**:150.

Neudeck *et al.* (2002) Expression variance, biochemical and immunological properties of *Toxoplasma gondii* dense granule protein GRA7. *Microbes Infect.* **4**:581.

Vercammen *et al.* (2000) DNA vaccination with genes encoding *Toxoplasma gondii* antigens GRA1, GRA7, and ROP2 induces partially protective immunity against lethal challenge in mice. *Infect. Immun.* **68**:38.

Ferguson *et al.* (1999) *In vivo* expression and distribution of dense granule protein 7 (GRA7) in the exoenteric (tachyzoite, bradyzoite) and enteric (coccidian) forms of *Toxoplasma gondii*. *Parasitology.* **119**:259.

Jacobs *et al.* (1999) Evaluation of recombinant dense granule antigen 7 (GRA7) of *Toxoplasma gondii* for detection of immunoglobulin G antibodies and analysis of a major antigenic domain. *Clin. Diagn. Lab. Immunol.* **6**:24.

Bonhomme *et al.* (1998) Quantitative immunolocalization of a P29 protein (GRA7), a new antigen of *toxoplasma gondii*. *J. Histochem. Cytochem.* **46**:1411.